

Product Number:	103591
Lot Number:	503753
Expiration:	28-Nov-2023

Component	Part Number	Lot Number
Bead Reagent	103581	233510
Detector Reagent	103583	233902
SBG Reagent	103585	233903
Control 1	103589	234101
Control 2	103590	234101
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	2.00 x 10 ⁷
Bead % Monomeric	88%
Isingle	150

Review/Approval

Sonia Khillar

QC

Sonia Khillar 19 DEC 2022

Name

Title

Signature/Date